

WAY-200070

Estrogen receptor ligand

Selective estrogen receptor ER β agonist displaying 68-fold greater selectivity over ER α (EC₅₀ = 2 and 155nM, respectively). It stimulates galanin expression in GnRH neurons and displays anxiolytic/antidepressant activity in various *in vivo* rodent models (administration to mice at 30mg/kg s.c.).

Citations: 4

[View Online »](#)

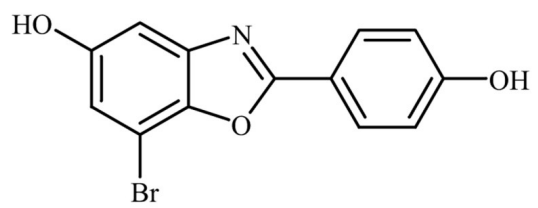
Ordering Information

[Order Online »](#)

BML-GR247-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	7-Bromo-2-(4-hydroxyphenyl)-1,3-benzoxazol-5-ol
Appearance	Off-white solid.
CAS	440122-66-7
Couple Target	Estrogen receptor
Couple Type	Ligand
Formula	$C_{13}H_8BrNO_3$
MW	306.1
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (12mg/ml) or 100% ethanol (2.5mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)